
PAK4

Product code: 07-126

Full-length human PAK4 [1-591(end) amino acids of accession number NP_005875.1] was expressed as N-terminal GST-fusion protein (91 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: SGKtide
ATP (μ M) Km app / Bin	: 2.5 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 12
IC50 at 1 mM ATP (nM)	: n.a.

